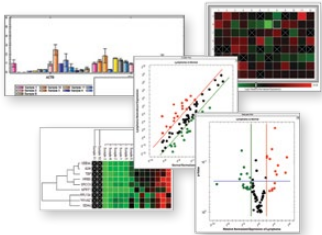




## Scalable System Maximizes Data Collection



The CFX Automation System II works with all CFX Real-Time PCR Detection Systems to enable walk-away, high-throughput quantitative PCR (qPCR) in a simplified format. This robotic plate handler for automated PCR facilitates the generation of large volumes of data and rapid data analysis.

### Features and Benefits of the CFX Automation System II

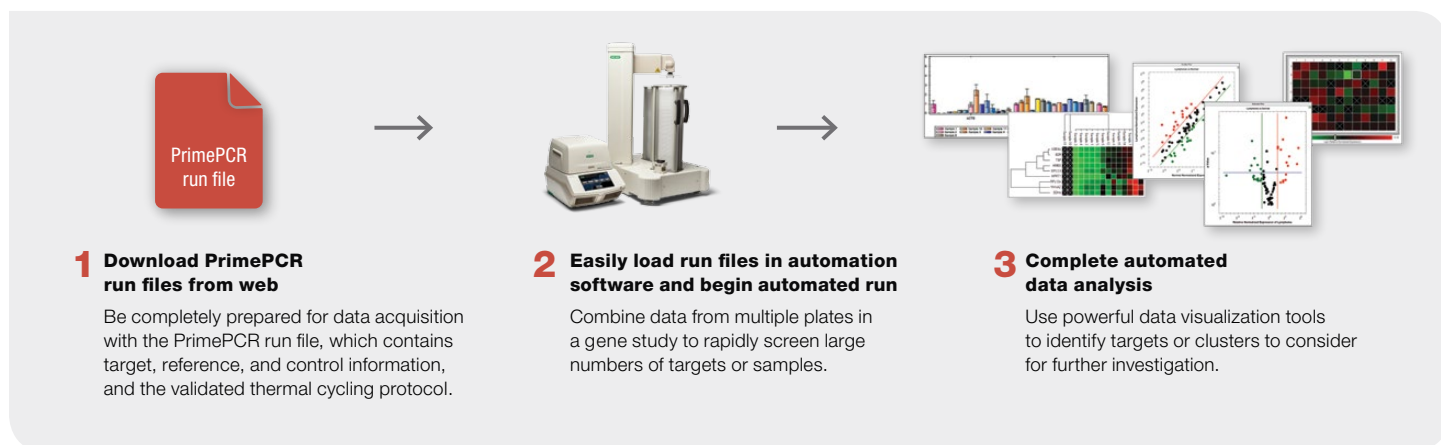
- **Save valuable laboratory space** — control 2 CFX Systems with a single plate handler in a compact footprint
- **Meet changing throughput demands** — begin with 1 real-time system and scale up to 2 systems; mix and match 96- and 384-well CFX Systems
- **Have confidence in your results** — use with PrimePCR™ Assays, which are expertly designed and wet-lab validated for proven performance; each assay for the human, mouse, and rat genomes was experimentally tested for optimal efficiency, specificity, sensitivity, and linear dynamic range
- **Analyze data when and where you want** — receive email notification with an attached data file at the completion of each run

Visit [bio-rad.com/web/CFXauto](http://bio-rad.com/web/CFXauto) for more information.



## Automating PrimePCR Is Simple with the CFX Automation System II

The automated PrimePCR workflow minimizes time-consuming setup, streamlines high-throughput data collection, and enables automated analysis of large data sets.



### Intuitive PCR Software Controls the CFX Automation System II

The easy-to-use automation software makes high-throughput real-time PCR simple and straightforward. The software is designed to work specifically with CFX Systems, eliminating the need for expertise with automation systems. Features include:

- **Customizable settings** — tailor automation runs to your specific needs
- **Powerful software features** — pair cycling information with specific barcoded plates and import barcode information from laboratory information management systems (LIMS)
- **Intuitive calibration wizard** — add CFX Systems at any point or recalibrate the automation system when required

### Specifications

Number of compatible real-time systems	2
Dimensions with single CFX System (W x D x H)	91 x 77 x 97 cm (36 x 30 x 38 in.)
Dimensions with dual configuration of CFX Systems (W x D x H)	99 x 77 x 97 cm (39 x 30 x 38 in.)
Number of racks	3
Number of 384-well plates per rack	48
Number of 96-well plates per rack	32
Automation software	CFX Automation Control Software 2.0 or higher
CFX Manager™ Software	Version 3.1.1621 or higher
Operating systems	Windows 7, Windows 8

### Ordering Information

Catalog #	Description
184-5075	<b>CFX Automation System II</b> , includes plate handler and barcode scanner, mounting plate, automation software; does not include CFX Real-Time PCR Detection System
185-5196	<b>CFX96 Touch™ Real-Time PCR Detection System</b> , includes C1000 Touch™ Thermal Cycler Chassis, CFX96™ Optical Reaction Module, CFX Manager Software, license for qbase+ Software, communication cable, reagents, consumables
185-5484	<b>CFX384 Touch™ Real-Time PCR Detection System</b> , includes C1000 Touch Thermal Cycler Chassis, CFX384™ Optical Reaction Module, CFX Manager Software, license for qbase+ Software, communication cable, reagents, consumables
HSP-9901	<b>Hard-Shell® Low-Profile 96-Well Skirted PCR Plates</b> , white shell, clear well, barcoded, 50
HSP-3905	<b>Hard-Shell 384-Well Standard PCR Plates</b> , clear shell, white well, barcoded, 50

Windows is a trademark of Microsoft Corporation.

Bio-Rad's real-time thermal cyclers are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 6,767,512 and 7,074,367.

Hard-Shell Plates are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 7,347,977; 6,340,589; and 6,528,302.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 3065 7550 **Canada** 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300 **Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026674 55 05 **Taiwan** 886 2 2578 7189 **Thailand** 1800 88 22 88 **United Kingdom** 020 8328 2000